

PRODUCT INFORMATION



tetranor-Misoprostol

Item No. 22693

Formal Name: 3-((1R,2R,3R)-3-hydroxy-2-((E)-4-hydroxy-4-methyloct-1-en-1-yl)-5-oxocyclopentyl)propanoic acid

MF: C₁₇H₂₈O₅

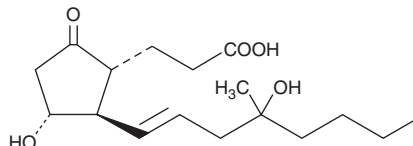
FW: 312.4

Purity: ≥95%

Supplied as: A solution in acetonitrile

Storage: -80°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

tetranor-Misoprostol is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of tetranor-misoprostol in ethanol and DMSO is approximately 50 mg/ml and approximately 100 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of tetranor-misoprostol is needed, it can be prepared by evaporating the acetonitrile and directly dissolving the neat oil in aqueous buffers. The solubility of tetranor-misoprostol in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

tetranor-Misoprostol is a metabolite of misoprostol (Item No. 13820).¹ It is formed when misoprostol undergoes rapid de-esterification to form misoprostol (free acid) (Item No. 13821), which is further metabolized to form the more polar tetranor-misoprostol in rats, dogs, and monkeys.

Reference

1. Schoenhard, G., Oppermann, J., and Kohn, F.E. Metabolism and pharmacokinetic studies of misoprostol. *Dig. Dis. Sci.* **30**(11 Suppl.), 126S-128S (1985).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/08/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM